


Product Number:	103208
Lot Number:	415915
Expiration:	7-Jun-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.33×10^9 Beads/mL

Review/Approval

ROBERT SOMOGYI QC MANAGER  23 AUG 2024

Name

Title

Signature/Date